

Zinquin ethyl ester (ultra pure)

Zinc dye

Zinquin ethyl ester is a lipophilic, zinc-sensitive, fluorescent derivative of Zinquin that is able to penetrate cell membranes. Cleavage of the ethyl ester group via cytosolic esterases in living cells impedes its efflux across the plasma membrane. It is selective for Zn (II) in the presence of physiological concentrations of Ca (II) and Mg (II) ions. Fluorescence of Zinquin ethyl ester is enhanced upon binding Zn (II), but the probe does not demonstrate a shift or change in emission wavelength. The probe is UV-excitable and emits in the blue region of the spectrum. Zinquin ethyl ester has been used to monitor intracellular zinc fluxes associated with apoptosis. Wavelength Maxima: Excitation 368nm, Emission 490nm

Citations: 4

[View Online »](#)

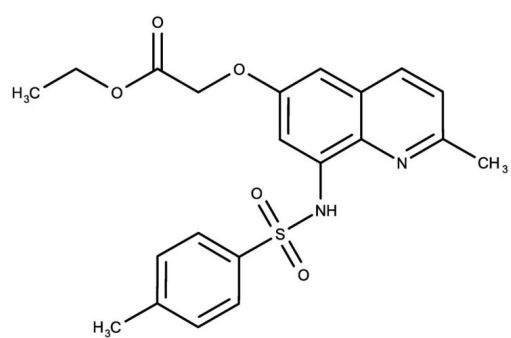
Ordering Information

[Order Online »](#)

ENZ-52151	5mg
-----------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Methyl-8-(4-methylphenylsulfonylamino)quinolinyl)oxyacetic acid ethyl ester
Appearance	White to pale-blue crystalline solid.
CAS	181530-09-6
Formula	$C_{21}H_{22}N_2O_5S$
MW	414.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO; slightly soluble in water.

Technical Info / Product Notes

This product is a member of the CELLESTIAL[®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL[®] reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com